

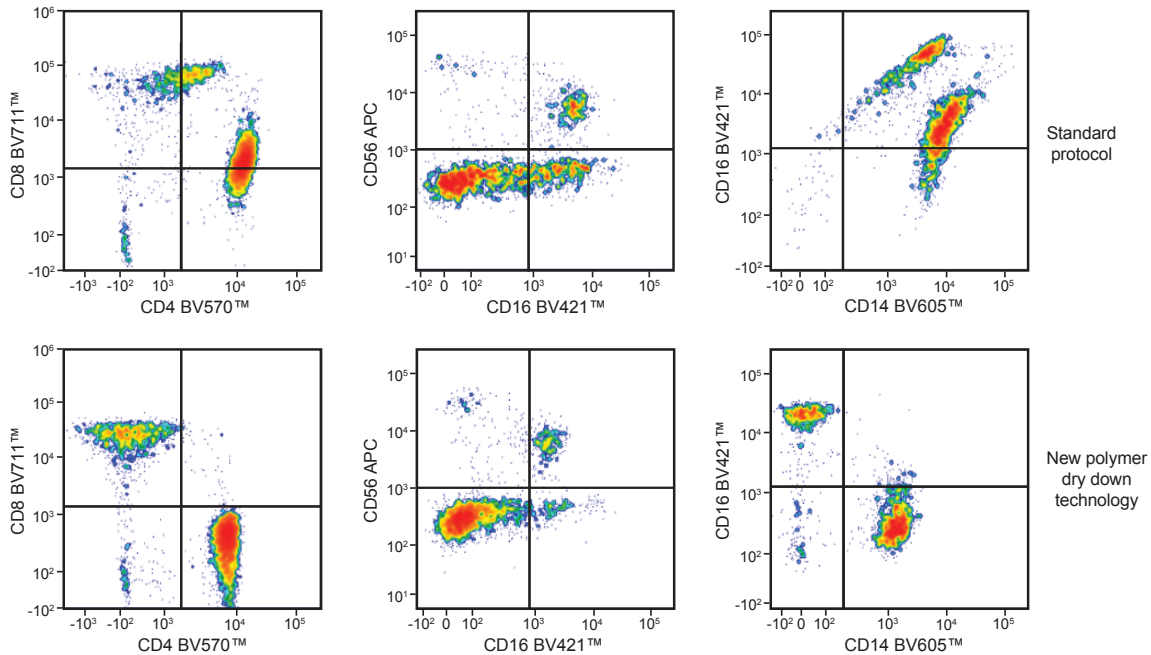
Multi-BV dry cocktails

Minimize polymer dye interactions

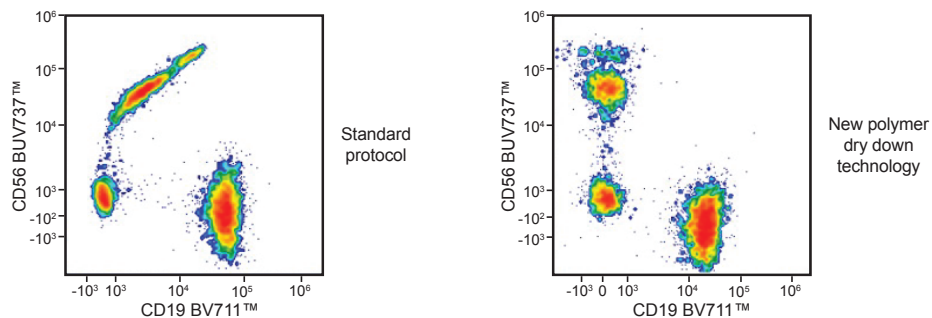
We now offer the option to include multiple Brilliant Violet™ dyes in our dry custom cocktails. These cocktails are formulated with our proprietary technology to minimize polymer dye interactions.

Learn more at [biolegend.com/en-us/custom-solutions](https://www.biolegend.com/en-us/custom-solutions)

Representative data



Human peripheral blood was stained with a 7-color antibody panel including BV421™, BV570™, BV605™, BV711™, BV750™, APC, and Spark Violet™ 538. Dot plots are gated on lymphocyte populations and depict dye-dye interaction (top plots), which are corrected in dry cocktails using BioLegend's new cocktail technology (bottom plots).



Human peripheral blood was stained with a 16-color antibody panel including seven Brilliant Violet™ dyes, BD Horizon™ BUV737, PE, PE/Cyanine7, APC/Fire™ 750, BD Horizon™ APC-R700, PE/Dazzle™ 594, and Alexa Fluor® 488. Dot plots are gated on lymphocyte populations and depict correction of spillover in dry cocktails using BioLegend's new cocktail technology.

Multi-BV dry cocktails

Flow cytometry cocktail comparison

	Liquid cocktails	Multi-BV dry cocktails
Product offering	Pre-formulated liquid cocktail	Pre-formulated dried cocktail
Capability	Conventional fluorophores plus one Brilliant Violet™ dye per vial	Conventional fluorophores, including multiple Brilliant Violet™ dyes
Accept non-BioLegend reagents	Yes	Yes
Custom conjugations	Yes	Yes
Stability	Varies based on cocktail composition	6–12 months
Storage conditions	2–8°C	20–25°C
Compensation tubes	Yes, free with evaluation material	Yes, free with evaluation material
FMO tubes	Yes, free with evaluation material	Yes, free with evaluation material
Evaluation tubes	Free of charge up to 20 tests; additional tubes may be requested at a small cost	Free of charge up to 20 tests; additional tubes may be requested at a small cost
QC data	Yes	Yes, pre- and post-dry down
Stimulation reagents spiked-in (PHA, LPS, etc.)	No	No
Packaging	Vials with custom sizes, labeling, and packing options	Custom packing formats available

We are unable to provide Brilliant Violet™ dyes to customers intending to use these reagents in a diagnostic or clinical setting.

Create your custom cocktail with us

Our goal is to optimize your workflow by increasing assay consistency with our custom cocktail reagents. Our dedicated lab and field scientists will work directly with you to customize reagents to your exact specifications to ensure outstanding quality and reproducibility.

Contact us to get started at [biolegend.com/en-us/custom-solutions](https://www.biolegend.com/en-us/custom-solutions)



BioLegend products are manufactured in an ISO 13485:2016 certified facility following GMP compliant procedures, ensuring the highest quality and stability.

BioLegend is a registered trademark of BioLegend, Inc. All other trademarks are the property of their respective owners.

BioLegend, Inc | www.biolegend.com | 8999 BioLegend Way | San Diego, CA 92121 | 1.877.246.5343 (877-BIOLEGEND) | International: 1.858.768.5800

07-0181-00